

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT CRF2 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1041

Lot Number: DC1041-072214

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human CRHR2 cDNA (GenBank Accession Number NM_001883.3) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

Background: CRF2 is a member of the corticotrophin-releasing factor (CRF) receptor family. It is expressed in the brain, blood vessels and intestine. CRF plays an integral role in the coordinating endocrine and behavioral responses to stress as well as in the pathophysiology of several neuropsychiatric diseases such as depression, anxiety and addiction. CRF2 is involved in stress responses, cardiovascular function and gastric motility. Recent clinical data suggest that CRF-related agents may be promising in the treatment of various endocrine, psychiatric, neurologic and inflammatory diseases. CRF2 agonists may be useful in the treatment of upper-GI inflammatory diseases.

Application: Functional assays

Figure 1

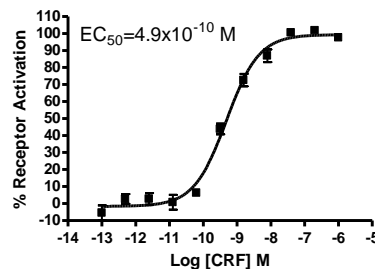


Figure 2

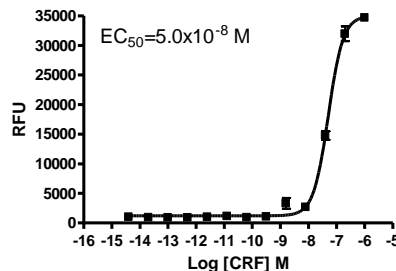


Figure 1. Dose-dependent inhibition of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). No transient transfection of Gα16.

References:

De Souza (1995) Corticotrophin-releasing factor receptors: physiology, pharmacology, biochemistry and role in central nervous system and immune disorders. *Psychoneuroendocrinology* 20:789-819.

Gravanis *et al.* (2005) The corticotrophin-releasing factor. *Curr Med Chem* 12:1503-1512.

Liaw *et al.* (1996) Cloning and characterization of the human corticotrophin-releasing factor-2 receptor complementary deoxyribonucleic acid. *Endocrinology* 137:72-77.

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.